

CUTANA™ Cloud

Upload. Analyze. Discover.



Get Started at EpiCypher.com/Cutana-Cloud

CUTANA™ Cloud is a web platform that **simplifies bioinformatic analysis of CUT&RUN and CUT&Tag data**. Designed for speed, security, and ease-of-use, it rapidly aligns raw sequencing files to quickly determine the success of an experiment.

Analysis of epigenomic data can be a major bottleneck, delaying experimental validation and slowing the path to biological insights. **CUTANA™ Cloud solves these challenges**, providing an accessible solution for every researcher — no coding required.

Why Choose CUTANA™ Cloud:

- **Fast & Reliable**
Results in as little as 90 minutes.
- **Biologist-Friendly**
No coding needed, intuitive UI.
- **High Security**
Enterprise-grade infrastructure.
- **Insight Ready**
Automated workflows.
- **NO Subscriptions**
Flexible, pay-as-you-go pricing.

CUTANA™ Cloud: Built for Every Researcher



Wet Lab Biologists

Steep learning curve for developing pipelines and accessing compute resources



Get results in as little as 90 minutes with just a few clicks



Bioinformaticians

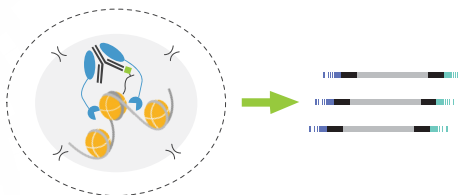
Routine analyses take time and pipelines are difficult to standardize



Standardized pipelines simplify workflows, with tools for custom analyses

Plug right into the CUTANA™ Ecosystem

DAY 1-3



CUTANA™ CUT&RUN /
CUT&Tag Assays

DAY 3



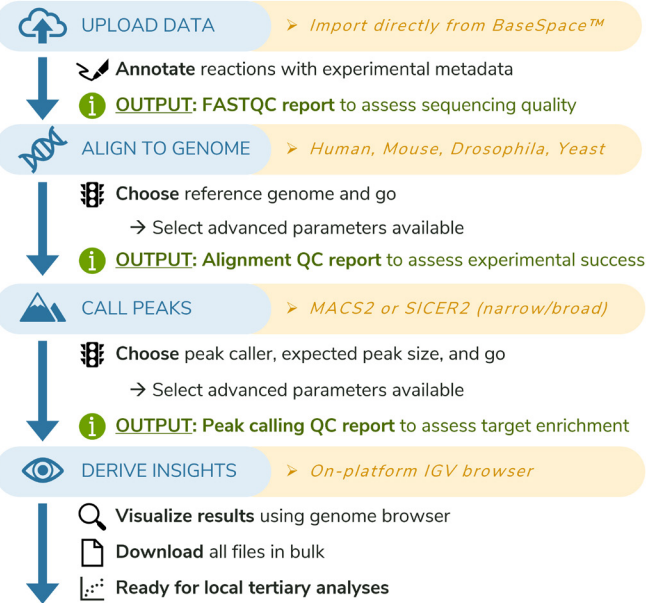
Next-Generation
Sequencing

DAY 4

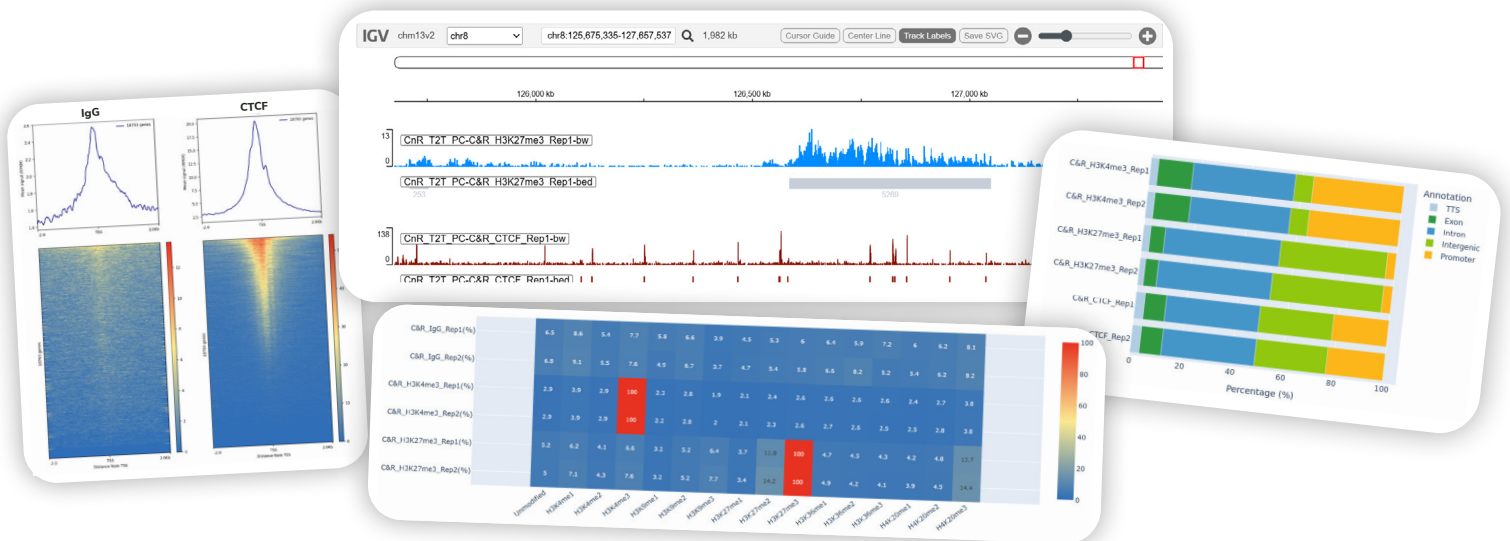


CUTANA™ Cloud
Data Analysis

The Cloud Workflow: From Data Upload to Discovery



Try CUTANA™ Cloud: Explore Sample Data for FREE



Start Fast — No contracts, No subscriptions.

Buy credits to analyze data at your own pace.

Each credit covers one genome alignment for a CUT&RUN or CUT&Tag reaction (up to 20 million reads). Each additional 20 million reads consumes an extra 0.25 credits. Peak calling is free.

Visit [EpiCypher.com/Cutana-Cloud](https://epicypher.com/Cutana-Cloud)

*For Research Use Only. Available to U.S. customers only.